

PRODUCT INFORMATION



Adipoyl-Coenzyme A (sodium salt)

Item No. 27858

Formal Name: S-(hydrogen hexanedioate)

coenzyme A, sodium salt

Synonyms: Adipoyl-CoA, Adipyl-CoA

MF: C₂₇H₄₄N₇O₁₉P₃S • XNa

FW: 895.7

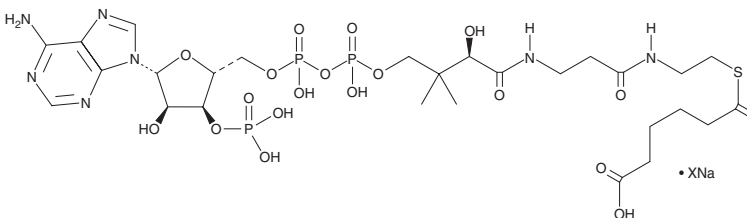
Purity: ≥70%

UV/Vis.: λ_{max}: 258 nm

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Adipoyl-coenzyme A (sodium salt) is supplied as a solid. Aqueous solutions of adipoyl-coenzyme A (sodium salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of adipoyl-coenzyme A (sodium salt) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Adipoyl-coenzyme A (adipoyl-CoA) is a medium-chain acyl CoA and an intermediate in the peroxisomal β -oxidation of dicarboxylic acids.¹ It is also an intermediate in the biosynthesis of adipic acid in genetically modified *E. coli*.²

References

1. Westin, M.A., Hunt, M.C., and Alexson, S.E. The identification of a succinyl-CoA thioesterase suggests a novel pathway for succinate production in peroxisomes. *J. Biol. Chem.* **280**(46), 38125-38132 (2005).
2. Yu, J.L., Xia, X.X., Zhong, J.J., et al. Direct biosynthesis of adipic acid from a synthetic pathway in recombinant *Escherichia coli*. *Biotechnol. Bioeng.* **111**(12), 2580-2586 (2014).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/17/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM